



Your Product

Quick to Tox

Quick to Clinic

GMP VHH Technology Platform

From cell bank to GMP in 8 months

With over **40 years of experience** in **recombinant protein** expression, we've developed a **robust microbial platform** dedicated to the **rapid and reliable production of VHH** antibodies in *E. coli* or *P. pastoris* (*K. phaffii*), with **scalable** output and full **GMP** compliance.



From RCB to clinic in
only **8 months**



End-to-End
Assistance



GMP Quality



Designed for
Tox & Ph I

VHH platform for optimized production

We've built a **robust VHH production platform**, developed through **extensive testing** on a diverse set of VHH models, including:

- | **Mono-, bi-, and tri-specific** formats
- | **Isoelectric points between pI 5 and 9**
- | **Diverse hydrophobicity** properties

This toolbox allows us to **rapidly assess your VHH product** against our **optimized fermentation and purification platform**, enabling the fast delivery of high-quality material suitable for **Tox studies** and **Phase I** human clinical trials.

The VHH toolbox

At the heart of our microbial platform

It integrates various **microbial host systems, fermentation strategies, purification**, and **QC methods** to rapidly fine-tune each production process.

Two expression platforms

- *P. pastoris* (*K. phaffii*)
- *E. coli*

Fermentation

- High-cell density fed-batch
- Control of critical fermentation parameters
- Animal-free media

Purification

- Ion-Exchange
- Hydrophobic Interaction Chromatography (HIC)
- Mix mode
- Affinity

Quality Control & Release

- GMP QC package
- Additional custom QC development available
- QP release

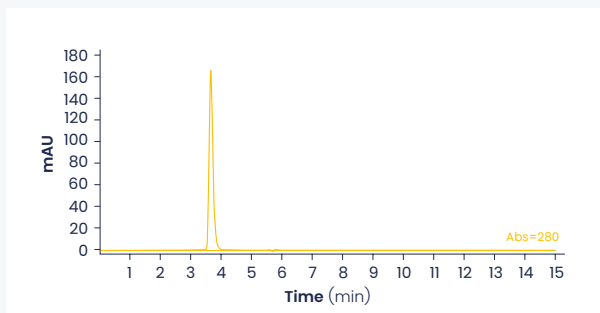
Support

- Project Management
- Dedicated team
- VHH Experts
- Tailored made solution



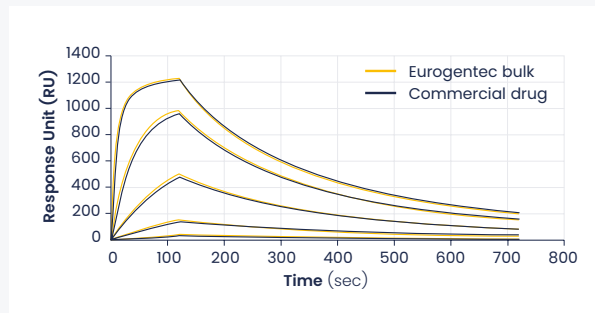
Attributes	Testing
Appearance	Colorimetry
	Turbidimetry
Physico-chemical	pH
Optional QCs	
Identity	ESI-Q-TOF
	RP-HPLC
Purity	RP-HPLC
Physico-chemical	Visible particles
	Osmolarity
Potency	Antigen binding assay by SPR
	ELISA

High quality



High Purity level

SE-HPLC analysis of a **trivalent bispecific VHH** shows a purity level of **98.7%**.



Binding affinity testing

SPR analysis (Biacore system) confirms that Eurogentec's bulk material displays **binding affinity equivalent to the commercial reference**, with K_d values of 19.7 nM and 20.9 nM, respectively.



Dedicated Support

We offer **personalized support**, **transparent communication**, and **tailored solutions** to meet your technical and regulatory needs. Our **dedicated project managers** and experienced **VHH specialists** guide you from early development to GMP release.

Contact us

For more information or to discuss your VHH project, feel free to reach out.



Scan and write us at
GMP-VHH@eurogentec.com